

15-Deoxy- Δ 12,14-prostaglandin J2

PPAR γ ligand

Prostaglandin D2 derived nuclear peroxisome proliferator-activated receptor γ (PPAR γ) ligand. Induces apoptosis in a number of cell types including human lung cancer cells, synoviocytes and macrophages. Promotes differentiation of fibroblasts to adipocytes. Antiinflammatory activity. Inhibits angiogenesis. Prevents TNF α induced leptin increase in mice. TRPA1 agonist. Serious doubts as to the biological relevance of this prostaglandin have been raised. Endogenous levels seem to be too low to have any effect as a PPAR γ ligand.

Citations: 18

[View Online »](#)

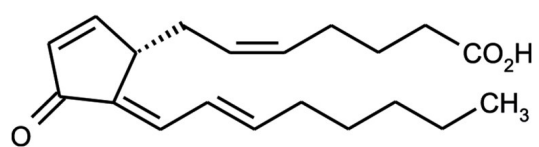
Ordering Information

[Order Online »](#)

BML-PG050-0001	1mg
BML-PG050-0005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for at least 1 year after receipt when stored, as supplied, at -20°C. Stock solutions are stable for up to 3 months at -20°C.
Handling	It is recommended to add an appropriate solvent directly to the vial.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	15-deoxy- Δ 12,14-PGJ2
Appearance	Neat oil. Vials may appear empty because the oil tends to coat the inside of the vial.
CAS	87893-55-8
Couple Target	PPAR, TRP channel
Couple Type	Ligand
Formula	$C_{20}H_{28}O_3$
MW	316.5
Purity	≥97% (TLC)
Solubility	Soluble in 100% ethanol (25mg/ml) or DMSO (25mg/ml).

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)